

Hyper Broth™

Hyper Broth™ is an APF Certified™ proprietary media formulation developed by AthenaES™ for improved expression of recombinant proteins in E. coli. The medium is a complex, rich formulation which supplies amino acids, vitamins, and glucose at high levels. This medium yields the highest level of biomass of all AthenaES™ media. The medium is buffered at pH 7.3 ± 0.2 and has the highest buffering capacity of Athena's Expression Media, which permits high glucose levels. For expression of recombinant proteins, induction should be done at cell densities three times that of LB Broth. Cell yields are typically 5 times that of LB Broth depending on the strain. Plasmid copy numbers are significantly higher than they are for LB Broth. This medium is suitable as the base for fermentation cultures and is excellent for feed-batch applications. Biomass yields will depend on the process and the strain employed. Hyper Broth™ powder is shipped with the necessary Glucose Nutrient Mix.

Ordering Information

[Order Online »](#)

AES-0107	500g
----------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Formulation Powder.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com